

# PRODUCT INFORMATION



## De-Boc-Cabazitaxel

Item No. 40116

**CAS Registry No.:** 1638286-64-2  
**Formal Name:** ( $\alpha$ R, $\beta$ S)- $\beta$ -amino- $\alpha$ -hydroxy-benzenepropanoic acid, (2aR,4S,4aS,6R,9S,11S,12S,12aR,12bS)-12b-(acetyloxy)-12-(benzoyloxy)-2a,3,4,4a,5,6,9,10,11,12,12a,12b-dodecahydro-11-hydroxy-4,6-dimethoxy-4a,8,13,13-tetramethyl-5-oxo-7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxet-9-yl ester

**Synonym:** Cabazitaxel Impurity IV

**MF:** C<sub>40</sub>H<sub>49</sub>NO<sub>12</sub>

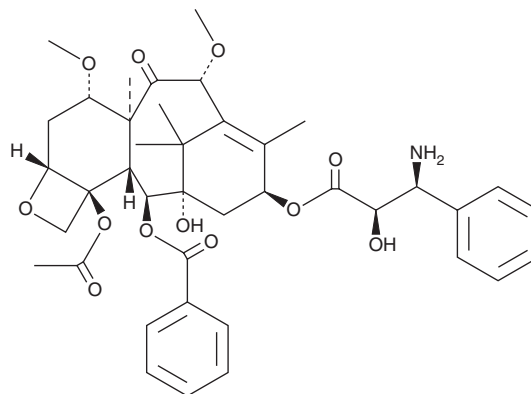
**FW:** 735.8

**Purity:**  $\geq$ 95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

De-Boc-Cabazitaxel is supplied as a solid. A stock solution may be made by dissolving the De-Boc-Cabazitaxel in the solvent of choice, which should be purged with an inert gas. De-Boc-Cabazitaxel is slightly soluble in methanol and DMSO.

### Description

De-Boc-Cabazitaxel is a potential impurity in commercial preparations of the semisynthetic taxane cabazitaxel.<sup>1</sup> It has also been used in the synthesis of tubulin-targeting fluorescent dyes.<sup>2</sup>

### References

1. Wang, Y., Feng, F., Chen, L., *et al.* Isolation, identification and characterization of potential impurities in cabazitaxel and their formation. *Magn. Reson. Chem.* **52(12)**, 783-788 (2014).
2. Grimm, F., Nizamov, S., and Belov, V.N. Green-emitting rhodamine dyes for vital labeling of cell organelles using STED super-resolution microscopy. *Chembiochem* **20(17)**, 2248-2254 (2019).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/07/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM